

## Product Information Sheet

### **Ordering Information**

Product Number: V100065

Product Name: Anti-Alpha-fetoprotein antibody \*Mouse anti-human, monoclonal IgG1\*

Unit Size: 50 ug

Storage Conditions: Freeze (<-15 °C)

Expiration Date: 12 months upon receiving

### **Chemical and Spectral Properties**

Appearance: Solid

Molecular Weight: N/A

Soluble In: Water

Excitation Wavelength: N/A

Emission Wavelength: N/A

### **Application Notes**

Alpha-fetoprotein (AFP) is a major plasma protein produced by the yolk sac and the liver during fetal development. It is widely used as diagnostic marker to screen for various developmental abnormalities and in adults, pancreatic cancer, lung cancer, liver cancer and liver cirrhosis. During pregnancy, high concentrations of AFP in the amniotic fluid may indicate severe congenital defects of a fetus, teratoma and fetal death. Anti-AFP monoclonal antibodies are highly specific for the detection of human AFP immunogens. They can be used for a broad range of immunoassays chemiluminescence immunoassays (CLIA) and rapid sandwich immunoassays.